

Product Number:	103400
Lot Number:	502921
Expiration:	16-Apr-2022
Platform(s):	SR-X

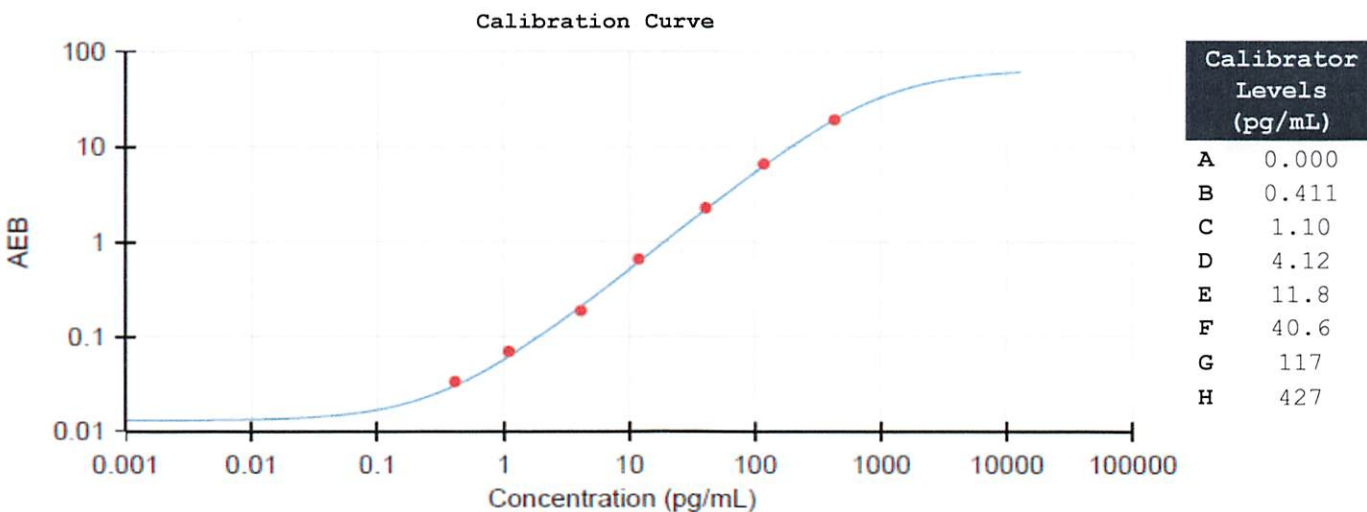


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive

Component	Part Number	Lot Number
Bead Reagent	102246	111804
Detector Reagent	102248	111715
SBG Reagent	102250	111406
Sample Diluent	103399	110609
Calibrators	102255	112301
Control 1	102256	112302
Control 2	102257	112302
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

Release Materials				Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	112302	4.57	2.28-5.08	
Control 2	112302	228	169-257	
Panel 1	008501	13.5	8.59-20.5	
Panel 2	008501	9.55	7.39-15.1	
Panel 3	008501	64.9	43.8-91.8	
Panel 4	009401	195	144-253	



Review/Approval _____

Junpil Park

QC Supervisor

 19 May 2021